

## General info

### Protocol information

Protocol name	Viral_NA_Duo
Modified by	admin
Kit name	MagJET Viral DNA and RNA Kit
Description	-

## Sample layout

1	
2	
3	
4	
5	
6	
7	
8	
9	
10	
11	
12	

## Reagent info







A		Elution strip		
Name	Well volume [µl]	Total reagent volume [µl]	Type	
Elution	100	-	Reagent	
A (Sample)		Viral NA plate		
Name	Well volume [µl]	Total reagent volume [µl]	Type	
Sample	200	-	Sample	
Lysis Buffer	200	-	Reagent	
Proteinase K	50	-	Reagent	
B (Tip)		Viral NA plate		
Name	Well volume [µl]	Total reagent volume [µl]	Type	
-	-	-	-	
C		Viral NA plate		
Name	Well volume [µl]	Total reagent volume [µl]	Type	
-	-	-	-	
D		Viral NA plate		
Name	Well volume [µl]	Total reagent volume [µl]	Type	
-	-	-	-	
E (Wash 1)		Viral NA plate		
Name	Well volume [µl]	Total reagent volume [µl]	Type	
Wash Buffer 1	700	-	Reagent	
F (Wash 2)		Viral NA plate		
Name	Well volume [µl]	Total reagent volume [µl]	Type	
Wash Buffer 2	800	-	Reagent	
G (Wash 3)		Viral NA plate		
Name	Well volume [µl]	Total reagent volume [µl]	Type	
Wash Buffer 2	800	-	Reagent	
H		Viral NA plate		
Name	Well volume [µl]	Total reagent volume [µl]	Type	
-	-	-	-	







## Dispensed reagents

A (Sample)		Viral NA plate		
Name	Step	Well volume [µl]	Total reagent volume [µl]	
Magnetic Beads	Pause	25	-	



## Steps data

 Tip	KingFisher Duo 12 tip comb		
 Pick-Up	Viral NA plate	(B) - Tip	
 Lysis	Viral NA plate	(A) - Sample	
Beginning of step	Precollect	No	
	Release beads	No	
Mixing / heating:	Mixing time, speed	00:15:00, Fast	
	Heating temperature [°C]	56	
End of step	Postmix	No	
	Collect beads	No	
	Post-temperature	No	
 Pause	Viral NA plate	(A) - Sample	
	Message	Add Magnetic Beads and 100% isopropanol	
Reagent(s)	Dispensing volume [µl]	475	
	Name	Magnetic Beads	
	Volume [µl]	25	
	Name	100% isopropanol	
	Volume [µl]	450	
 Bind	Viral NA plate	(A) - Sample	
Beginning of step	Precollect	No	
	Release beads	Yes	
Mixing / heating:	Shake 1 time, speed	00:00:30, Bottom mix	
	Shake 2 time, speed	00:00:30, Fast	
	Loop count	5	
	Heating during mixing	No	
End of step	Postmix	No	
	Collect count	3	
	Collect time [s]	3	
	Post-temperature	No	
 Wash 1	Viral NA plate	(E) - Wash 1	
Beginning of step	Precollect	No	
	Release time, speed	00:00:05, Fast	
Mixing / heating:	Mixing time, speed	00:02:00, Half mix	
	Heating during mixing	No	
End of step	Postmix	No	
	Collect count	3	
	Collect time [s]	2	
	Post-temperature	No	

	Wash 2	Viral NA plate	(F) - Wash 2
	Beginning of step	Precollect	No
		Release time, speed	00:00:05, Fast
	Mixing / heating:	Mixing time, speed	00:02:00, Half mix
		Heating during mixing	No
	End of step	Postmix	No
		Collect count	3
		Collect time [s]	2
		Post-temperature	No
	Wash 3	Viral NA plate	(G) - Wash 3
	Beginning of step	Precollect	No
		Release beads	No
	Mixing / heating:	Shake 1 time, speed	00:01:30, Slow
		Shake 2 time, speed	00:00:15, Bottom mix
		Loop count	5
		Heating during mixing	No
	End of step	Postmix	No
		Collect count	3
		Collect time [s]	1
		Post-temperature	No
	Dry	Viral NA plate	(G) - Wash 3
		Dry time	00:03:00
		Tip position	Outside well / tube
	Elution	Elution strip	(A)
	Beginning of step	Precollect	No
		Release time, speed	00:00:05, Fast
	Mixing / heating:	Shake 1 time, speed	00:05:00, Medium
		Shake 2 time, speed	00:02:00, Paused
		Tip position when paused	Tip edge in liquid
		Heating temperature [°C]	56
	End of step	Postmix	No
		Collect count	3
		Collect time [s]	2
		Post-temperature	No
	Release Beads	Viral NA plate	(G) - Wash 3
		Release time, speed	00:00:05, Fast
	Leave	Viral NA plate	(A) - Sample

## Lot info

No lot numbers have been defined.